

Product Number:	100376
Lot Number:	523314
Expiration:	21-Feb-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.35 - 7.45	PASS

RESULT

pH Reading	7.36
------------	------

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature]
29 Sep 2025
Signature/Date